

Kinase Assay Buffer II

Catalog # K02-09

Lot # L2027-6

Product Description

Stock solution of buffer used for assaying kinase protein.

Formulation

Frozen solution of buffer containing: 25 mM MOPS, pH7.2, 12.5mM β -glycerol-phosphate, 25mM $MgCl_2$, 12.5mM $MnCl_2$, and 2mM EDTA.

Add to final concentration of 0.25mM DTT (Catalog# D86-09) prior to use. DTT Solution supplied in separate vial.

Sterility

Buffer passed through a 0.2 μ filter membrane and packaged aseptically.

Storage and Stability

Store product at $-20^{\circ}C$. For optimal storage, aliquot buffer into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

Kinase Assay Buffer II

Catalog #	K02-09
Lot #	L2027-6
Stability	1 yr at $-20^{\circ}C$ from date of shipment
Storage & Shipping	Store product at $-20^{\circ}C$. For optimal storage, aliquot buffer into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped frozen.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.